

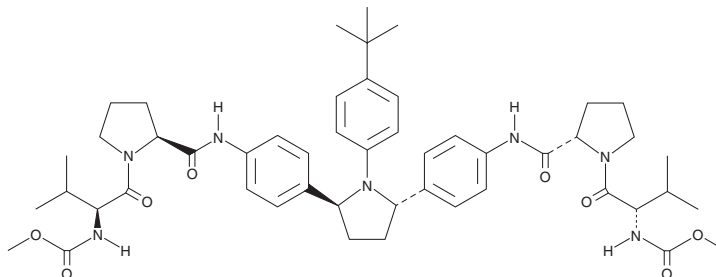
PRODUCT INFORMATION



Ombitasvir

Item No. 24116

CAS Registry No.: 1258226-87-7
Formal Name: 2,2'-[[[(2S,5S)-1-[4-(1,1-dimethylethyl)phenyl]-2,5-pyrrolidinediyl]di-4,1-phenylene]bis[N-(methoxycarbonyl)-L-valyl-L-prolinamide]
Synonym: ABT-267
MF: C₅₀H₆₇N₇O₈
FW: 894.1
Purity: ≥98%
UV/Vis.: λ_{max}: 252 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Ombitasvir is supplied as a crystalline solid. A stock solution may be made by dissolving the ombitasvir in the solvent of choice, which should be purged with an inert gas. Ombitasvir is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of ombitasvir in these solvents is approximately 30 mg/ml.

Ombitasvir is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ombitasvir should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Ombitasvir has a solubility of approximately 0.3 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Ombitasvir is an orally bioavailable and potent inhibitor of the hepatitis C virus (HCV) non-structural protein 5A (NS5A).^{1,2} It reduces HCV replication in stable replicon cell lines with EC₅₀ values ranging from 0.82 to 19.3 pM against genotypes 1a, 1b, 2a, 2b, 3a, 4a, and 5a and an EC₅₀ value of 366 pM for 6a.^{1,2} Against subgenomic replicons from clinical isolates of genotypes 1a-6a expressed in Huh-7 derived cells, ombitasvir inhibits HCV replication with EC₅₀ values ranging from 0.1 pM for genotype 4a to 68 pM for genotype 6a.² Formulations containing ombitasvir have been used in the treatment of HCV genotypes 1 and 4.

References

1. Zhang, X. Direct anti-HCV agents. *Acta Pharmaceutica Sinica B* **6**(1), 26-31 (2015).
2. Krishnan, P., Beyer, J., Mistry, N., et al. *In vitro* and *in vivo* antiviral activity and resistance profile of ombitasvir, an inhibitor of hepatitis C virus NS5A. *Antimicrob. Agents Chemother.* **59**(2), 979-987 (2015).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/31/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM